Glucagon monoclonal antibody (08)

Glucagon is a peptide hormone of 29 amino-acid residues produced by the alpha cells of the pancreatic islets. Its effect is to raise blood sugar as a stress hormone and counter-regulatory hormone for insulin. It is produced by cell-type-specific processing of proglucagon, which in the alpha cells is split to form glicentin-related pancreatic polypeptide, glucagon and major proglucagon fragment, while the same precursor is split to form glicentin, oxyntomodulin (C-terminally extended glucagon-37), GLP-1 and GLP-2 in the L cells of the intestinal mucosa.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BPD-ABS-058-08-1	1mg
BPD-ABS-058-08-04	400µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA

Clone 08

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M

sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Synthetic N-terminal fragment corresponding to human

glucagon coupled to carrier.

Isotype IgG1ĸ

Purity Detail Protein A-affinity purified.

Species Reactivity Human, Porcine, Rat

UniProt ID P01275

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

_

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com